Rose-β-D-Gal

Rose- β -D-Gal (6-Chloro-3-indoxyl- β -D-galactopyranoside) is a chromogenic substrate for β -galactosidase. The product is similar to X-gal, but generates an intense pink colored precipitate ($\lambda_{max} \sim 540$ nm) on enzymatic hydrolysis.



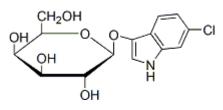
Product attributes

Call us: 800-304-5357 Email: btinfo@biotium.com

Product Description

Rose- β -D-Gal (6-Chloro-3-indoxyl- β -D-galactopyranoside) is a chromogenic substrate for β -galactosidase. The product is similar to X-gal, but generates an intense pink colored precipitate ($\lambda_{max} \sim 540$ nm) on enzymatic hydrolysis.

- White solid soluble in DMSO
- Store at 4°C and protect from light
- MW: 329.74
- C₁₄H₁₆CINO₆



References

- 1. Molecular Cloning: A laboratory Manual, Cold Spring Harbor Laboratory, B. 14, p186 (1989).
- 2. Biotechniques 27, 438 (1999).
- 3. Oncogene 10, 2323 (1995).
- 4. Dev Biol 161, 77 (1994).
- 5. J Histochem Cytochem 42, 1299 (1994).
- 6. BioTechniques 15, 974 (1993).
- 7. BioTechniques, 15, 292 (1993).

This datasheet was generated on November 4, 2025 at 02:04:49 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/rose-b-d-gal-6-chloro-3-indoxyl-b-d-galactopyranoside/